

PRODUCT INFORMATION



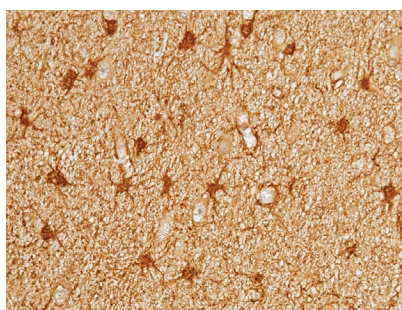
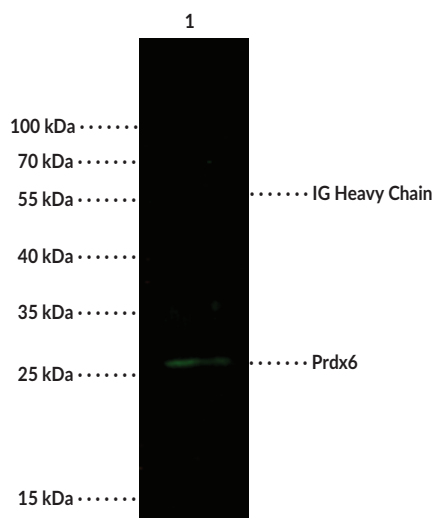
Peroxiredoxin-6 Monoclonal Antibody (Clone 03)

Item No. 43212

Overview and Properties

Contents:	This vial contains 100 µl of protein A-purified monoclonal antibody.
Synonyms:	aiPLA ₂ , Acidic Calcium-independent Phospholipase A ₂ , 1-Cys Peroxiredoxin, Glutathione-dependent Peroxiredoxin, LPCAT5, Non-selenium Glutathione Peroxidase, NSGPx, Prx6
Immunogen:	Recombinant human Prx6
Cross Reactivity:	(+) Prx6, (-) <i>E. coli</i> cell lysate
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-80°C (as supplied)
Stability:	≥1 year
Storage Buffer:	0.2 µm filtered solution in PBS, with 10% trehalose, pH 7.0
Clone:	03
Host:	Mouse
Isotype:	IgG1
Applications:	ELISA, Immunohistochemistry-paraffin (IHC-P), Immunoprecipitation (IP); the recommended starting dilution is 1:1,000-1:2,000 for ELISA and 1:50-1:200 for IHC-P; the recommended starting concentration is 4-6 µl/mg of lysate for IP. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Immunohistochemical staining of formalin-fixed and paraffin-embedded human Prdx6 in human brain using Peroxiredoxin-6 Monoclonal Antibody at a dilution of 1:60.

Lane 1: 0.5 mg HeLa whole cell lysate, 2 µl Peroxiredoxin-6 Monoclonal Antibody, and 15 µl of 50% Protein G agarose.

Prdx6 was immunoprecipitated using Peroxiredoxin-6 Monoclonal Antibody at a dilution of 1:500 followed by Dylight 800-labeled antibody to Mouse IgG (H+L) at a dilution of 1:7,500. It has an observed band size of 26 kDa under reducing conditions.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/02/2025

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Peroxiredoxin-6 (Prx6) is a 1-Cys-containing Prx that is involved in the maintenance of cellular thiol redox homeostasis and phospholipid metabolism.¹ It exists as a homodimer and is composed of a thioredoxin fold-containing catalytic domain, which contains a peroxidatic cysteine (Cp), and a C-terminal arm.² Unlike 2-Cys-containing Prxs, Prx6 does not contain a resolving cysteine (Cr) and uses glutathione (GSH) instead. Prx6 is ubiquitously expressed and localizes to the cytosol, nucleus, mitochondria, and plasma membrane.^{1,3} It has three enzymatic activities: glutathione peroxidase (GPX), acidic calcium-independent phospholipase A₂ (aiPLA₂), and lysophosphatidylcholine acyltransferase (LPCAT).^{1,4,5} The LPCAT activity, which selectively uses palmitoyl-CoA as a substrate, is coupled to the aiPLA₂ activity and is selective for lysophosphatidylcholine over lysophosphatidylethanolamine, lysophosphatidylglycerol, lysophosphatidylinositol, and lysophosphatidylserine.⁵ These activities, coupled with the GPX activity, provide Prx6 the ability to repair oxidized cell membranes.¹ Prx6 is also involved in the activation of NADPH oxidase 2 (NOX2). Tumor levels of Prx6 are increased in patients with non-small cell lung cancer (NSCLC) and serum levels of Prx6 are decreased in patients with asthma.⁶ Cayman's Peroxiredoxin-6 Monoclonal Antibody (Clone O3) can be used for ELISA, immunohistochemistry-paraffin (IHC-P), and immunoprecipitation (IP) applications.

References

1. Arevalo, J.A. and Vázquez-Medina, J.P. The role of peroxiredoxin 6 in cell signaling. *Antioxidants (Basel)* **7(12)**, 172 (2018).
2. Choi, H.J., Kang, S.W., Yang, C.H., *et al.* Crystal structure of a novel human peroxidase enzyme at 2.0 Å resolution. *Nat. Struct. Biol.* **5(5)**, 400-406 (1998).
3. Dammeyer, P. and Arnér, E.S. Human Protein Atlas of redox systems - What can be learnt? *Biochim. Biophys. Acta* **1810(1)**, 111-138 (2011).
4. Chen, J.W., Dodia, C., Feinstein, S.I., *et al.* 1-Cys peroxiredoxin, a bifunctional enzyme with glutathione peroxidase and phospholipase A₂ activities. *J. Biol. Chem.* **275(37)**, 28421-28427 (2000).
5. Fisher, A.B., Dodia, C., Sorokina, E.M., *et al.* A novel lysophosphatidylcholine acyl transferase activity is expressed by peroxiredoxin 6. *J. Lipid Res.* **57(4)**, 587-596 (2016).
6. Jia, W., Dong, C., and Li, B. Anti-oxidant and pro-oxidant effects of peroxiredoxin 6: A potential target in respiratory diseases. *Cells* **12(1)**, 181 (2023).